

PRODUCT INFORMATION



1,2-Stearoyl-3-trimethylammoniumpropane (chloride)

Item No. 37404

CAS Registry No.: 220609-41-6
Formal Name: N,N,N-trimethyl-2,3-bis[(1-oxooctadecyl)oxy]-1-propanaminium, monochloride

Synonyms: DSTAP, 18:0 TAP

MF: C₄₂H₈₄NO₄ • Cl

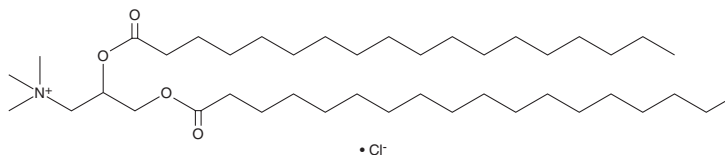
FW: 702.6

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

1,2-Stearoyl-3-trimethylammoniumpropane (DSTAP) (chloride) is supplied as a solid. A stock solution may be made by dissolving the DSTAP (chloride) in the solvent of choice, which should be purged with an inert gas. DSTAP (chloride) is slightly soluble in methanol.

Description

DSTAP is a cationic lipid that has been used in the formation of lipid nanoparticles (LNPs).¹ Intraperitoneal administration of LNPs containing DSTAP and encapsulating resiquimod (R-848; Item No. 14806) increases peritoneal fluid levels of IFN- α , the number of tumor-associated CD8⁺ T cells, and reduces peritoneal tumor luminescence in a CT26-luc colorectal cancer mouse model of colorectal cancer with peritoneal metastases.

Reference

1. Pauli, G., Chao, P.-H., Qin, Z., *et al.* Liposomal resiquimod for enhanced immunotherapy of peritoneal metastases of colorectal cancer. *Pharmaceutics* **13**(10), 1696 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/29/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM